

## Product Information

### High Sensitivity Glucose Assay Kit

Catalog Number **MAK181**

Storage Temperature  $-20^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Glucose is a primary energy source that naturally occurs in its free state in fruits and other plant parts. Abnormal glucose levels have been associated with several metabolic dysfunctions such as hypoglycemia, hyperglycemia, and diabetes mellitus. Measurements of glucose levels in tissues and body fluids (such as blood and urine) are often used for the diagnosis of glucose-related disorders. Glucose levels are also monitored to check the efficacy of therapeutics such as insulin and sulphonylureas in type 2 diabetics.<sup>1,2</sup>

This kit is a highly sensitive assay for determining glucose levels (ranging from 20–100 pmole/well) in a variety of biological samples. Glucose concentration is determined by a coupled enzyme assay, which results in a fluorometric ( $\lambda_{\text{ex}} = 535/\lambda_{\text{em}} = 587 \text{ nm}$ ) product proportional to the glucose present.

### Components

The kit is sufficient for 100 assays in 96 well plates.

Glucose Assay Buffer Catalog Number MAK181A	25 mL
Glucose Probe, in DMSO Catalog Number MAK181B	0.4 mL
Glucose Enzyme Mix Catalog Number MAK181C	1 vL
Glucose Substrate Mix Catalog Number MAK181D	1 vL
Glucose Standard, 100 mM Catalog Number MAK181E	0.1 mL

### Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate. It is recommended to use white plates with clear bottoms for fluorescence assays.
- Fluorescence multiwell plate reader.
- 10 kDa Molecular Weight Cut-Off (MWCO) Spin Filter (optional for protein-containing samples)

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

Glucose Assay Buffer – Allow buffer to come to room temperature before use.

Glucose Probe – Warm to room temperature prior to use to melt DMSO. Mix well by pipetting, then aliquot and store, protected from light and moisture, at  $-20^{\circ}\text{C}$ .

Glucose Enzyme Mix – Reconstitute in 220  $\mu\text{L}$  of Glucose Assay Buffer. Mix well by pipetting, then aliquot and store at  $-20^{\circ}\text{C}$ . Keep on ice while in use. Use within 2 months of reconstitution.

Glucose Substrate Mix – Reconstitute in 220  $\mu\text{L}$  of water. Mix well by pipetting, then aliquot and store at  $-20^{\circ}\text{C}$ . Keep on ice while in use. Use within 2 months of reconstitution.

Glucose Standard – Ready to use. Store at  $-20^{\circ}\text{C}$ .

### Storage/Stability

The kit is shipped on wet ice and storage at  $-20^{\circ}\text{C}$ , protected from light, is recommended.

### Procedure

All samples and standards should be run in duplicate.

#### Glucose Standards for Fluorometric Detection

Dilute 10  $\mu\text{L}$  of the 100 mM Glucose Standard Solution with 990  $\mu\text{L}$  of water to prepare a 1 mM (1 nmole/ $\mu\text{L}$ ) Standard Solution. Dilute 10  $\mu\text{L}$  of the 1 mM (1 nmole/ $\mu\text{L}$ ) Standard Solution into 990  $\mu\text{L}$  of water to prepare a 10  $\mu\text{M}$  (10 pmole/ $\mu\text{L}$ ) Standard Solution. Add 0, 2, 4, 6, 8 and 10  $\mu\text{L}$  of the 10  $\mu\text{M}$  Standard Solution into a 96 well plate, generating 0 (blank), 20, 40, 60, 80, and 100 pmole/well standards. Add Glucose Assay Buffer to each well to bring the volume to 50  $\mu\text{L}$ .

#### Sample Preparation

Liquid samples can be assayed directly.

Tissue samples (10 mg) or cells ( $1 \times 10^6$ ) should be homogenized on ice with 100  $\mu\text{L}$  of ice cold Glucose Assay Buffer. Centrifuge at 12,000 rpm for 5 minutes to remove insoluble material and collect the supernatant.

Proteins in samples may impact background levels. To remove proteins from samples, deproteinize using a 10 kDa MWCO spin filter.

For unknown samples, it is suggested to test several sample volumes to make sure the readings are within the standard curve range.

Add 1–50  $\mu\text{L}$  (1–10  $\mu\text{g}$ ) of samples into duplicate wells of a 96 well plate. Bring samples to a final volume of 50  $\mu\text{L}$  with Glucose Assay Buffer.

High levels of NADH can result in sample background. To correct for the background, include a Sample Blank for each sample by omitting the Glucose Enzyme Mix. The Sample Blank readings can then be subtracted from the sample readings.

### Assay Reaction

1. Set up the appropriate Reaction Mixes according to the scheme in Table 1. 50  $\mu\text{L}$  of Reaction Mix is required for each reaction (well).

**Table 1.**  
Reaction Mixes

Reagent	Samples and Standards	Sample Blank
Glucose Assay Buffer	45 $\mu\text{L}$	47 $\mu\text{L}$
Glucose Probe	1 $\mu\text{L}$	1 $\mu\text{L}$
Glucose Enzyme Mix	2 $\mu\text{L}$	–
Glucose Substrate Mix	2 $\mu\text{L}$	2 $\mu\text{L}$

2. Add 50  $\mu\text{L}$  of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting, and incubate the reaction for 30 minutes at  $37^{\circ}\text{C}$ . Protect the plate from light during the incubation.
3. Measure fluorescence intensity ( $\lambda_{\text{ex}} = 535/\lambda_{\text{em}} = 587 \text{ nm}$ ).

## Results

### Calculations

The background for the assays is the value obtained for the 0 (blank) Glucose Standard. Correct for the background by subtracting the blank standard value from all readings. Background values can be significant and must be subtracted from all readings. Use the values obtained from the appropriate Glucose Standards to plot a standard curve.

Note: A new standard curve must be set up each time the assay is run.

Subtract the Sample Blank value from the sample reading to obtain the corrected fluorescence measurement. Using the corrected fluorescence measurement, determine the amount of glucose present in the sample from the standard curve.

### Concentration of Glucose

$$S_a/S_v = C$$

where:

$S_a$  = Amount of Glucose in sample well (pmole) from standard curve

$S_v$  = Sample volume ( $\mu$ L) added to well

$C$  = Concentration of Glucose in sample

Glucose molecular weight: 180.16 g/mole

### Sample Calculation

Amount of Glucose ( $S_a$ ) = 25.84 pmole  
(from standard curve)

Sample volume ( $S_v$ ) = 50  $\mu$ L

Concentration of Glucose in sample:

$$25.84 \text{ pmole}/50 \text{ } \mu\text{L} = 0.5168 \text{ pmole}/\mu\text{L}$$

$$0.5168 \text{ pmole}/\mu\text{L} \times 180.16 \text{ pg}/\text{pmole} = 93.11 \text{ pg}/\mu\text{L}$$

### References

1. Golden, S.H., and Sapir, T., Methods for insulin delivery and glucose monitoring in diabetes: summary of a comparative effectiveness review. J. Manag. Care Pharm., **18(6 suppl.)**, S1–17 (2012).
2. Intensive blood-glucose control with sulphonylureas or insulin compared with conventional treatment and risk of complications in patients with type 2 diabetes (UKPDS 33). UK Prospective Diabetes Study (UKPDS) Group. Lancet, **352(9131)**, 837–853 (1998).

**Troubleshooting Guide**

<b>Problem</b>	<b>Possible Cause</b>	<b>Suggested Solution</b>
Assay not working	Cold assay buffer	Assay Buffer must be at room temperature
	Omission of step in procedure	Refer and follow Technical Bulletin precisely
	Plate reader at incorrect wavelength	Check filter settings of instrument
	Type of 96 well plate used	For fluorescence assays, use white plates with clear bottoms.
Samples with erratic readings	Samples prepared in different buffer	Use the Assay Buffer provided or refer to Technical Bulletin for instructions
	Samples were not deproteinized	Use 10 kDa MWCO Spin Filters to deproteinize samples
	Cell/Tissue culture samples were incompletely homogenized	Repeat the sample homogenization, increasing the length and extent of homogenization step.
	Samples used after multiple freeze-thaw cycles	Aliquot and freeze samples if samples will be used multiple times
	Presence of interfering substance in the sample	If possible, dilute sample further
	Use of old or inappropriately stored samples	Use fresh samples and store correctly until use
Lower/higher readings in samples and standards	Improperly thawed components	Thaw all components completely and mix gently before use
	Use of expired kit or improperly stored reagents	Check the expiration date and store the components appropriately
	Allowing the reagents to sit for extended times on ice	Prepare fresh Reaction Mixes before each use
	Incorrect incubation times or temperatures	Refer to Technical Bulletin and verify correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly
Non-linear standard curve	Use of partially thawed components	Thaw and resuspend all components before preparing the Reaction Mixes
	Pipetting errors in preparation of standards	Avoid pipetting small volumes
	Pipetting errors in the Reaction Mix	Prepare Reaction Mixes whenever possible
	Air bubbles formed in well	Pipette gently against the wall of the plate well
	Standard stock is at incorrect concentration	Refer to the standard dilution instructions in the Technical Bulletin
	Calculation errors	Recheck calculations after referring to Technical Bulletin
	Substituting reagents from older kits/lots	Use fresh components from the same kit
Unanticipated results	Samples measured at incorrect wavelength	Check the equipment and filter settings
	Samples contain interfering substances	If possible, dilute sample further
	Sample readings above/below the linear range	Concentrate or dilute samples so readings are in the linear range

VI,KVG,LS,MAM 09/14–1